

Mononucleosome, (H3.1ΔN32), 167x601 DNA, Recombinant Human

Catalog No. 16-2130
Lot No. 19252001
Pack Size 50 µg



EpiCypher®

Product Description:

Mononucleosomes assembled from recombinant histones expressed in *E. coli* (two each of histones H2A, H2B, H3 and H4. Accession numbers: H2A-P04908; H2B-O60814; H3.1-P68431; H4-P62805), with the amino acid sequence of H3 beginning with glycine 33 (amino acids 1-32 are deleted), wrapped in a 167bp DNA sequence containing a 147bp Lowary and Widom 601 positioning sequence. The 601 sequence is flanked by a 10 bp sequence. The nucleosome is the basic subunit of chromatin.

Formulation:

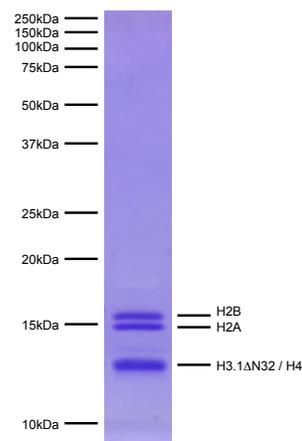
Purified recombinant mononucleosomes (50 µg total mass, 25.1 µg protein and 24.9 µg DNA) in 59 µl 10 mM Tris-HCl pH 7.5, 1 mM EDTA, 25 mM NaCl, 2 mM DTT and 20% glycerol. Concentration of nucleosomes is 4.09 µM. Nucleosome molecular weight = 207,312.13 Da.

Storage and Stability:

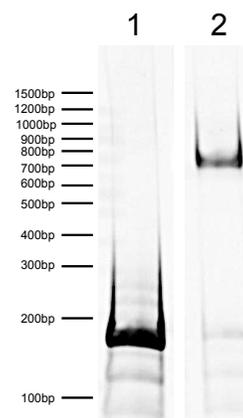
Stable for six (6) months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

Mononucleosomes, (H3.1ΔN32), Recombinant Human are highly purified and are suitable for use as substrates in enzyme screening assays or for nucleosome binding experiments. The absence of post-translational histone modifications makes them ideal for conducting enzyme activity and screening assays. **EpiCypher Mononucleosomes, Recombinant do not contain free DNA which could alter assayed activities.**



Protein Gel Data: Coomassie stained PAGE gel of proteins in Mononucleosome, (H3.1ΔN32), 167x601, Recombinant Human (1 µg) demonstrates the purity of the histones in the preparation. Sizes of molecular weight markers and positions of the core histones (H2A, H2B, H3.1ΔN32, and H4) are indicated. H3.1ΔN32 and H4 co-migrate.



DNA Gel Data: Mononucleosome, (H3.1ΔN32), 167x601 DNA, Recombinant Human resolved via native PAGE and stained with ethidium bromide to visualize DNA. **Lane 1:** Free DNA (100 ng). **Lane 2:** Intact nucleosomes (400 ng).

This product is for *in vitro* research use only and is not intended for use in humans or animals.